



August 25, 2021

From: Eric Giesing, Director

Re: LS 13 320 Instrument Obsolescence

Dear LS 13 320 Customer,

This is an obsolescence notice for the LS 13 320 particle size analyzer. Due to a combination of parts obsolescence and a strategic business pivot, the instrument will become obsolete and no longer available for purchase as of August 25, 2021. We will honor and ship instruments for purchase orders received up to the obsolescence date of August 25, 2021.

There is not currently, nor is there projected to be, a replacement or alternative product available from Beckman Coulter Life Sciences. Please note that after the date of obsolescence, additional software and hardware improvements or updates will not be made.

We will continue to fully service and support the LS 13 320 for a period of five (5) years after the obsolescence date up to August 25, 2026. During this time, calibrations, repairs, and reagents required for routine service will be available, however, breakdown repairs cannot be guaranteed if the required repair component is made obsolete by our supplier.

Instrument part numbers affected by the LS 13 320 obsolescence include:

6605638	AQUEOUS LIQUIDS MODULE
383497	KIT, UPGRADE, SONICATION, ULM2
383665	UNIVERSAL LIQUIDS MODULE 2
6605211	LS SAMPLE CELL PIDS
6605530R	VAR SPEED MODULE PLUS, REFURB
6605686	LS AUTO-PREP STATION, (NC)
7801054	R-CELL ASY. MICRO VOLUME
7809200	LS, VACUUM CLEANER (OPTION)
8310540	KIT, ACCESSORY LS13320
8310542	KIT, PACKAGING, DPS/ULM
8310573	KIT, ACCESSORY, ULM
8310693	KIT, V CHECK, LS13 320
A86579	PRGMD CD ROM, LS 3 SERIES, V6.03, ENABLED, CNTRL SWRE
A86580	PRGMD CD ROM, LS 3 SERIES, .01, CNTRL SWRE
A88098	DATA DISPLAY, 20.1 INCH TFT ACTIVE MATRIX LCD DISPLAY W/ USB HUB
B00262	SALES GROUP TN, COMPUTER AND MONITOR FOR LS AND MS SERIES

Instrument accessory part numbers affected by the LS 13 320 obsolescence include:

383497	KIT, UPGRADE, SONICATION, ULM2
383665	UNIVERSAL LIQUIDS MODULE 2
6605211	LS SAMPLE CELL PIDS
6605530R	VAR SPEED MODULE PLUS, REFURB
6605686	LS AUTO-PREP STATION, (NC)
7801054	R-CELL ASY. MICRO VOLUME
7809200	LS, VACUUM CLEANER (OPTION)
8310540	KIT, ACCESSORY LS13320
8310542	KIT, PACKAGING, DPS/ULM
8310573	KIT, ACCESSORY, ULM
8310693	KIT, V CHECK, LS13 320
A86579	PRGMD CD ROM, LS 3 SERIES, V6.03, ENABLED, CNTRL SWRE
A86580	PRGMD CD ROM, LS 3 SERIES, V6.01, CNTRL SWRE
A88098	DATA DISPLAY, 20.1 INCH TFT ACTIVE MATRIX LCD DISPLAY W/ USB HUB
B00262	SALES GROUP TN, COMPUTER AND MONITOR FOR LS AND MS SERIES

We apologize for the abrupt nature of this decision and are happy to help answer questions you may have moving forward. Please contact your local sales representative to find out how this may impact a current or planned order. For technical questions or details regarding continued service and support, please contact Customer Technical Support (CTS) at (800) 369-0333.

Thank you,



Eric Giesing
 Director, Biotechnology Business Unit, Beckman Coulter Life Sciences